

Product Number:	101736 (103159)
Lot Number:	501653 (90380)
Expiration:	7-Feb-2020
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler

Associate Director, QC

P. Stadler

12 Feb 2019

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.